

# Complex I Human Protein Quantity Dipstick Assay Kit ab109722

[8 References](#) [3 Images](#)

### Overview

<b>Product name</b>	Complex I Human Protein Quantity Dipstick Assay Kit
<b>Sample type</b>	Cell culture extracts, Tissue
<b>Assay type</b>	Sandwich (quantitative)
<b>Species reactivity</b>	<b>Reacts with:</b> Cow, Human
<b>Product overview</b>	<p>ab109722 contains 48 dipsticks and necessary components to quantify the levels of the fully assembled Complex I enzyme complex in human and bovine samples. The kit includes sufficient materials to generate a standard curve and evaluate several unknown samples.</p> <p>Based on the immunologic sandwich assay, the kit utilizes two monoclonal antibodies (mAbs) specific to different antigens present on the Complex I enzyme complex. One antibody is immobilized on the nitrocellulose membrane in a thin line perpendicular to the length of the dipstick, while the other is gold-conjugated and combined with the sample mix. The sample contents containing the gold-conjugated mAb wick past the mAb immobilized on the dipstick. When assembled Complex I is present in the sample, a red line appears at the site of the anti-Complex I antibody line. The signal intensity is directly related to the amount of Complex I in the sample. The signal intensity is best measured by a dipstick reader or may be analyzed by another imaging system. To identify defects in Complex I that do not affect enzyme assembly combine this assay with <a href="#">ab109720</a> to determine the relative specific activity of Complex I.</p>
<b>Notes</b>	<p>All components are stable in their provided containers at room temperature out of direct sunlight.</p> <p>After diluting the 10X Blocking Buffer to 2X, store at 4°C. For long-term storage, all buffers can be stored at 4°C.</p>
<b>Tested applications</b>	<b>Suitable for:</b> Sandwich ELISA
<b>Platform</b>	Reagents

### Properties

**Storage instructions** Store at +4°C. Please refer to protocols.

Components	48 tests
Buffer B (10X Blocking solution)	1 x 0.4ml

Components	48 tests
Dipsticks	1 x 30 units
Extraction Buffer (ab260490)	1 x 15ml
Gold-conjugated antibody (dried in microplate wells)	1 x 30 tests
Wash buffer	1 x 2ml

## Applications

**The Abpromise guarantee** Our **Abpromise guarantee** covers the use of ab109722 in the following tested applications. The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
<b>Sandwich ELISA</b>		Use at an assay dependent dilution.

## Images

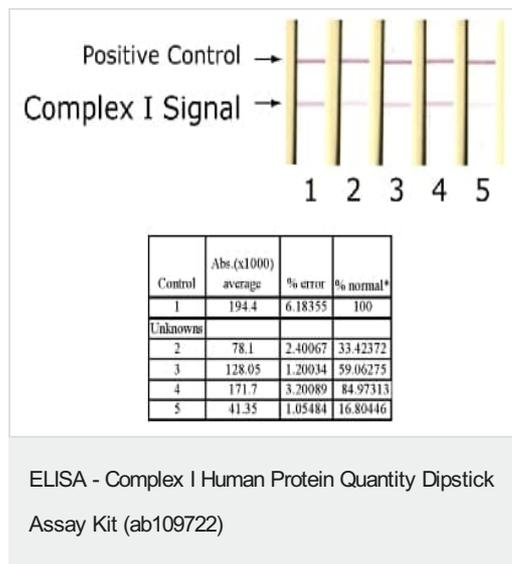


Figure 2. An example using ab109722 to quantify Complex I levels using various concentration of human fibroblast extract.

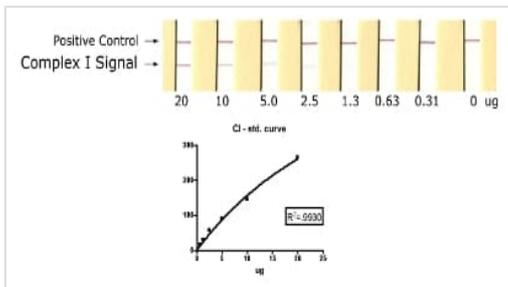
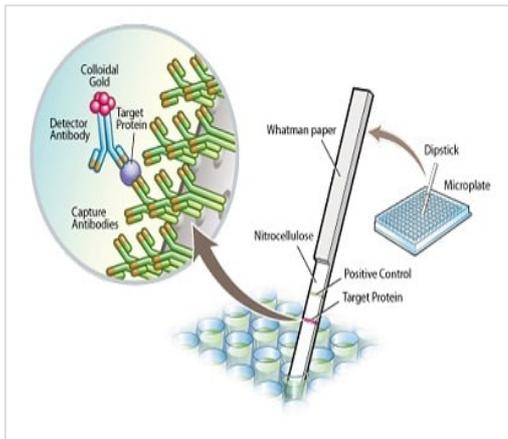


Figure 1. An example using ab109722 to quantify Complex I levels using various concentration of human fibroblast extract.

ELISA - Complex I Human Protein Quantity Dipstick Assay Kit (ab109722)



Sandwich ELISA - Complex I Human Protein Quantity Dipstick Assay Kit (ab109722)

Dipstick assays use the well-established lateral flow concept, whereby capture antibodies are striped onto nitrocellulose membrane and a Whatman paper wicking pad draws the sample through the antibody bands. Detector antibodies, conjugated to gold, are dried in the wells of a 96-well plate. Sample is added to the well, the dipstick inserted, and within minutes the line for each target is revealed as the protein-detector antibody-gold complex binds with the capture antibodies. Multiplexing dipstick assays have multiple target protein lines. A positive control goat anti-mouse antibody line is included on all assays to ensure that adequate wicking of the sample occurred.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors